

Diphtheria toxin

(*Corynebacterium diphtheriae*)

Elongation factor 2 (EF-2) catalyzes the GTP hydrolysis-dependent translocation reaction that is responsible for the movement of the ribosome along mRNA during protein synthesis. Diphtheria toxin catalyzed ADP-ribosylation of EF-2 inhibits the translocation reaction which halts protein synthesis and ultimately results in cell death. Diphtheria toxin is cell permeable and can translocate across endosomal membranes in response to low pH3.

Citations: 10

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-G135-0001	1mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°C
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Formulation	Lyophilized.
Reconstitution	Reconstitute with 0.5mL of sterile distilled water
Source	From <i>Corynebacterium diphtheriae</i> .

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)